
CHCHD2 Polyclonal Antibody

(Catalog # A73604)

Background

The protein encoded by this gene belongs to a class of eukaryotic CX(9)C proteins characterized by four cysteine residues spaced ten amino acids apart from one another. These residues form disulfide linkages that define a CHCH fold. In response to stress, the protein translocates from the mitochondrial intermembrane space to the nucleus where it binds to a highly conserved 13 nucleotide oxygen responsive element in the promoter of cytochrome oxidase 4I2, a subunit of the terminal enzyme of the electron transport chain. In concert with recombination signal sequence-binding protein J, binding of this protein activates the oxygen responsive element at four percent oxygen. In addition, it has been shown that this protein is a negative regulator of mitochondria-mediated apoptosis. In response to apoptotic stimuli, mitochondrial levels of this protein decrease, allowing BCL2-associated X protein to oligomerize and activate the caspase cascade. Pseudogenes of this gene are found on multiple chromosomes. Alternative splicing results in multiple transcript variants.

Description

CHCHD2 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

Q9Y6H1

Purification

Affinity Purification

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 75-145 of human CHCHD2 (NP_057223.1).

Storage

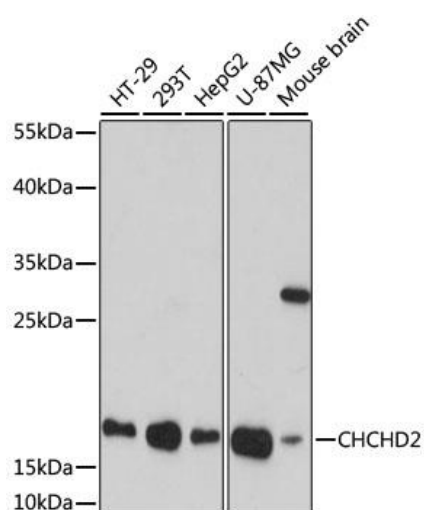
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

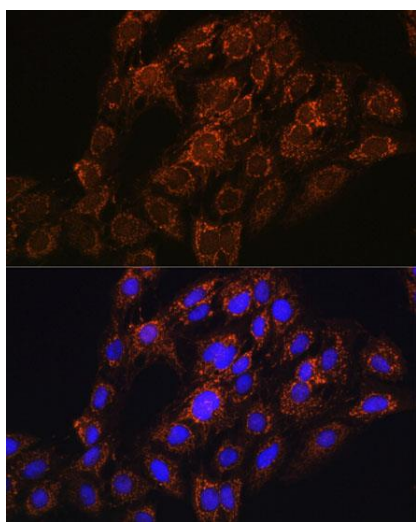
CHCHD2; C7orf17; MNRR1; NS2TP; PARK22

Application

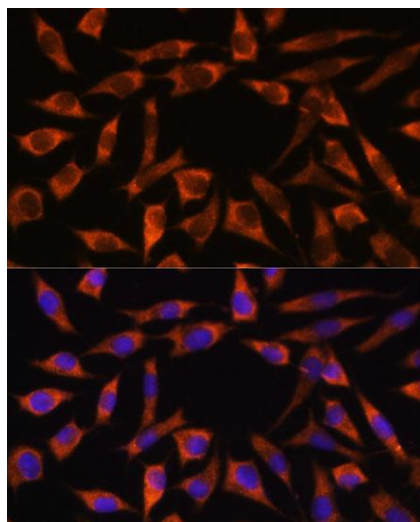
WB, IHC, IF; Recommended dilution: WB: 1:500-1:2000, IHC: 1:50-1:200, IF:1:50-1:200



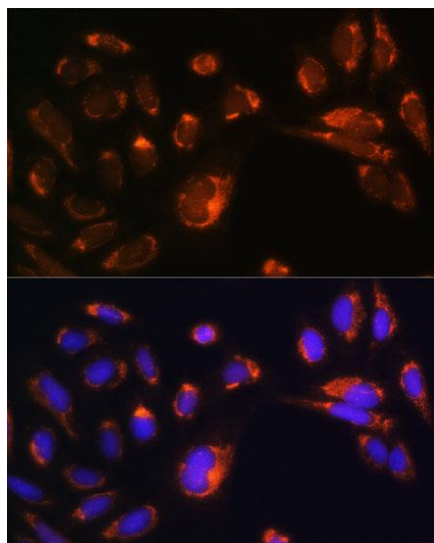
Western blot analysis of extracts of various cell lines, using CHCHD2 Polyclonal Antibody at 1:1000 dilution. Secondary Antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution. Lysates/proteins: 25ug per lane. Blocking buffer: 3% nonfat dry milk in TBST. Exposure time: 90s.



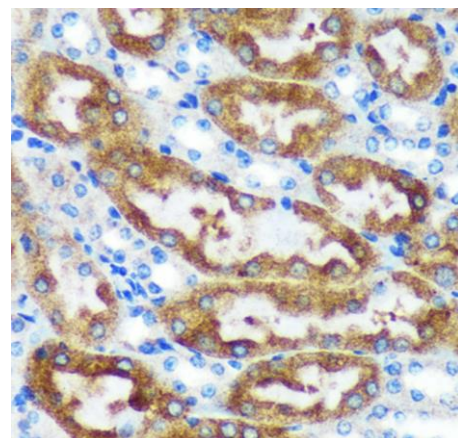
Immunofluorescence analysis of C6 cells using CHCHD2 Polyclonal Antibody at dilution of 1:100. Blue: DAPI for nuclear staining.



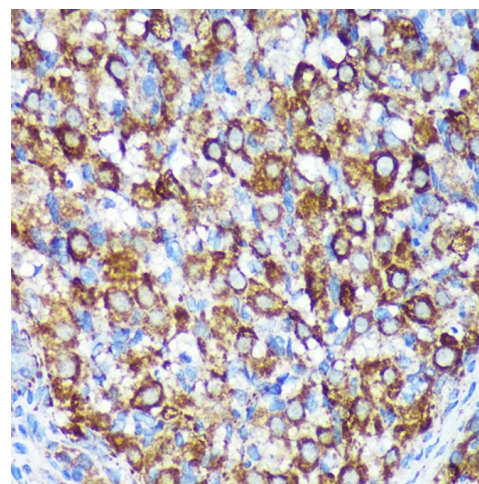
Immunofluorescence analysis of L929 cells using CHCHD2 Polyclonal Antibody at dilution of 1:100. Blue: DAPI for nuclear staining.



Immunofluorescence analysis of U-2 OS cells using CHCHD2 Polyclonal Antibody at dilution of 1:100. Blue: DAPI for nuclear staining.



Immunohistochemistry of paraffinembedded Mouse kidney using CHCHD2 Polyclonal Antibody at dilution of 1:100 (40x lens).



Immunohistochemistry of paraffinembedded Rat ovary using CHCHD2 Polyclonal Antibody at dilution of 1:100 (40x lens).